

---

## Ubiquitin Monoclonal Antibody [P4D1]

[Catalog No. A-8415-100]

---

**Description**

Mouse monoclonal antibody to Ubiquitin, Clone P4D1

**Size**

100 µg

**Formulation**

Liquid. In PBS containing 0.01% sodium azide.

**Immunogen**

Denatured bovine ubiquitin.

**Specificity**

Species Independent

**Isotype**

IgG1

**Purification**

Ion exchange purified

**Storage**

Store at -20°C. Avoid multiple freeze/thaw cycles.

**Application**

WB (1:1000)

*This product is for research purposes only. Not intended for use in diagnostic procedures.*